

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
MOUSE RECOMBINANT GPR119 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DCm1298-1

Lot Number: DCm1298-1-011917

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Cellbanker 2

Host cell: CHO-K1

Transfection: Expression vector containing full-length mouse GPR119 cDNA (GenBank Accession Number: BC115558) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM-F12, 10% FBS

Stability: 1 – 2 days after thawing

Background: GPR119 is a G protein-coupled receptor for the endogenous lipid signaling agent oleoylethanolamide (OEA). GPR119 is predominantly expressed in the pancreas (beta-cells) and gastrointestinal tract (enteroendocrine cells). In vitro studies have indicated a role for GPR119 in the modulation of insulin release by pancreatic beta-cells and of GLP-1 secretion by gut enteroendocrine cells.

Application: Functional assays

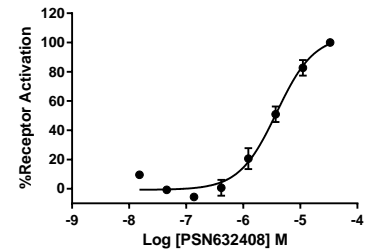
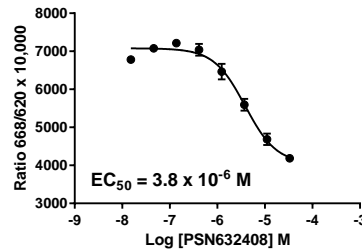


Figure Legend: Dose-dependent accumulation of intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Hilary *et al.* (2006) Deorphanization of a G protein-coupled receptor for oleoylethanolamide and its use in the discovery of smallmolecule hypophagic agents. *Cell Metabolism* 3:167-175.

Overton *et al.* (2007) GPR 119, a novel G protein-coupled receptor target for the treatment of type 2 diabetes and obesity. *Br J Pharmacol* 153 suppl 1:S76-S81.

FOR RESEARCH USE ONLY.

All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.